

Product Number:	103526
Lot Number:	217812
Expiration:	27-Jun-2023
Storage:	2 - 8°C

Laboratory Analysis

Characteristic	Means of Analysis	Result
Concentration	Multisizer	1.42 x 10 ⁹ Beads/mL

Review/Approval

Sonia Khillan
Name

QC Manager
Title

Sonia Khillan / 10 OCT 2022
Signature/Date